

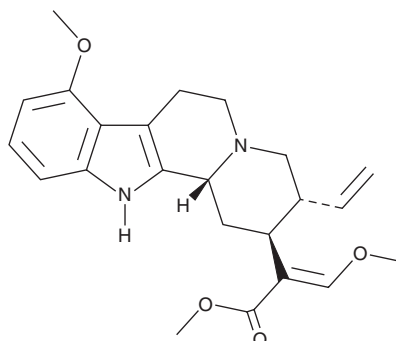
PRODUCT INFORMATION



Isopaynantheine

Item No. 39879

CAS Registry No.: 22032-51-5
Formal Name: (αE,2S,3R,12bR)-3-ethenyl-1,2,3,4,6,7,12,12b-octahydro-8-methoxy-α-(methoxymethylene)-indolo[2,3-a]quinolizine-2-acetic acid, methyl ester
Synonym: 3-Isopaynantheine
MF: C₂₃H₂₈N₂O₄
FW: 396.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Isopaynantheine (Item No. 39879) is an analytical reference standard categorized as an opioid.¹ Isopaynantheine is an alkaloid found in *M. speciosa* (Kratom in Thai).² It has antinociceptive activity in mice.¹ This product is intended for research and forensic applications.

References

1. Chakraborty, S., Uprety, R., Daibani, A.E., *et al.* Kratom alkaloids as probes for opioid receptor function: Pharmacological characterization of minor indole and oxindole alkaloids from kratom. *ACS Chem. Neurosci.* **12**(14), 2661-2678 (2021).
2. Philipp, A.A., Wissenbach, D.K., Weber, A.A., *et al.* Metabolism studies of the Kratom alkaloids mitraciliatine and isopaynantheine, diastereomers of the main alkaloids mitragynine and paynantheine, in rat and human urine using liquid chromatography- linear ion trap- mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **879**(15-16), 1049-1055 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/24/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM